

**Quantifoil™ R 2/2 on 200 copper mesh**

Cat. No.	Amount
X-103-Cu200	100 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a

**Applications:**

Support film for biological samples in Cryo-EM techniques

**Description:**

QUANTIFOIL™ R 2/2 carbon film has circular holes of 2 µm diameter with an interspace of 2 µm in both dimensions, applied on 200 copper mesh. Suitable for magnifications between 30.000x and 40.000x.